

ZytoLight® SPEC

ALK/EML4 TriCheck™

Probe

For the detection of ALK-EML4 rearrangements by fluorescence *in situ* hybridization (FISH).

The ZytoLight SPEC ALK/EML4 TriCheck™ Probe is designed to detect inversions involving the chromosomal region 2p23.1-p23.2 harboring the ALK gene and the chromosomal region 2p21 harboring the EML4 gene. Moreover, using this probe it is possible to discriminate between EML4-ALK inversions and translocations affecting ALK, but not EML4.

Ordering Information

[Order Online »](#)

ZTV-Z-2117-50	50µl
ZTV-Z-2117-200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application FISH

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), using 10 µl probe solution per test.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).
For the Manual please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com